DMNPE-Caged D-Luciferin

DMNPE-caged D-luciferin (D-Luciferin, 1-(4,5-dimethoxy-2-nitrophenyl)ethyl ester) is a D-luciferin ester derivative which can cross cell membranes efficiently. 1,2 Once inside the cells, the ester is continuously hydrolyzed to a supply of



Product attributes

Call us: 800-304-5357

CAS number

223920-67-0

Product Description

D-luciferin.

DMNPE-caged D-luciferin (D-Luciferin, 1-(4,5-dimethoxy-2-nitrophenyl)ethyl ester) is a D-luciferin ester derivative which can cross cell membranes efficiently. 1.2 Once inside the cells, the ester is continuously hydrolyzed to a supply of D-luciferin. Alternatively a burst of D-luciferin is generated by UV photolysis.

- Light yellow solid soluble in DMSO
- Store at -20°C and protect from light
- C₂₁H₁₉N₃O₇S₂
- MW: 489.52

References

- 1. Lett Appl Microbiol 30, 223 (2000).
- 2. Biotechniques 15, 848 (1993).

This datasheet was generated on December 3, 2025 at 01:06:04 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dmnpe-caged-d-luciferin-d-luciferin-1-45-dimethoxy-2-nitrophenylethyl-ester/